

PRODUCT INFORMATION



PF-9363

Item No. 37778

CAS Registry No.: 2569009-58-9

Formal Name: 2,6-dimethoxy-N-[4-methoxy-6-(1H-pyrazol-1-ylmethyl)-1,2-benzisoxazol-3-yl]-benzenesulfonamide

Synonym: CTx-648

MF: $C_{20}H_{20}N_4O_6S$

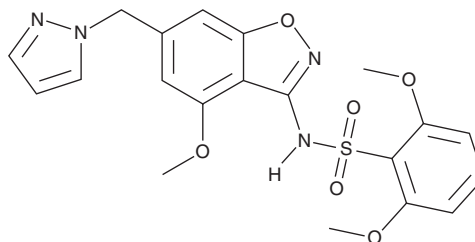
FW: 444.5

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PF-9363 is supplied as a solid. A stock solution may be made by dissolving the PF-9363 in the solvent of choice, which should be purged with an inert gas. PF-9363 is soluble in the organic solvent DMSO.

Description

PF-9363 is an inhibitor of lysine acetyltransferase 6A (KAT6A), also known as monocytic leukemia zinc finger protein (MOZ) and KAT6B, also known as MOZ-related factor (MORF; K_i s = 0.27 and 2.4 nM, respectively).¹ It is selective for KAT6A and KAT6B over KAT7, KAT5, and KAT8 (K_i s = 70, 420, and 670 nM, respectively). PF-9363 reduces the proliferation of, and the levels of acetylation of histone 3 lysine 23 (H3K23Ac) and lysine 14 (H3K14Ac) in, ZR-75-1 breast epithelial carcinoma cells (IC_{50} s = 0.37, 0.85, and 120 nM, respectively). It reduces tumor levels of H3K23Ac and tumor volume without affecting body weight in a ZR-75-1 breast cancer mouse xenograft model when administered at doses of 0.03, 0.3, or 5 mg/kg per day. PF-9363 (1 mg/kg per day) reduces tumor volume and increases survival in a patient-derived xenograft (PDX) mouse model of breast cancer.

Reference

1. Sharma, S., Chung, C.-Y., Uryu, S., *et al.* Discovery of a highly potent, selective, orally bioavailable inhibitor of KAT6A/B histone acetyltransferases with efficacy against KAT6A-high ER+ breast cancer. *Cell Chem. Biol.* **30** (2023).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/12/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM